

## Data Sheet

### CDK11A/CyclinD3, GST-tags

Human, Recombinant, N-terminal GST-tags

**Catalog #:** 100595

**Lot #:** 191216

**Conc.:** 0.1 mg/ml

**Formulated in:** 50 mM Tris-HCl, pH 7.5, 50 mM NaCl, 10 mM glutathione, 0.25 mM DTT, 0.1 mM EDTA, 0.1 mM PMSF, 25% glycerol.

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles.

**Description:** CDK11A/Cyclin D3 complex of human CDK11A, also known as CDC2L2, p58GTA, and PITSLRE, GenBank Accession No. NM\_024011, a.a. 361-502(end) with N-terminal GST-tag, MW= 73 kDa, and human Cyclin D3, also known as CCND3, GenBank Accession No. NM\_001760, a.a. 1-292(end) with N-terminal GST-tag, MW= 58 kDa. Co-expressed in a Baculovirus infected Sf9 cell expression system.

#### References:

1. Lahti, J. M. *et al*, *Nature Genet.* **7**: 370-375, 1994.
2. Bunnell, B. *et al*, *Proc. Nat. Acad. Sci.* **87**: 7467-7471, 199

## Quality Assurance

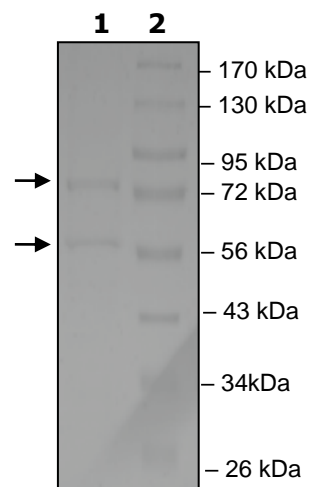
### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
CDK11A/Cyclin D3

**Lane 2:**  
Protein Marker

**Purity:**  $\geq 80\%$

**MW:**  
CDK11A: 73 kDa  
CyclinD3: 58 kDa



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)